

Product Number:	101473
Lot Number:	401614
Expiration:	16-Jul-2025
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.38	7.35 - 7.45	PASS

Review/Approval

ROBERT SOMOGYI  
Name

QC MANAGER  
Title

 19 JAN 2024  
Signature/Date